

ADP-Gel / ADP- Agarose

ADP immobilized on agarose gel

Adenosine-5'-O-diphosphate (ADP) has been immobilized on a polymeric matrix via the terminal phosphate. May be suitable for the affinity purification of ADP-binding proteins. BIOLOG also offers immobilized cytidine-5'-O-diphosphate (CDP-Gel, Cat. No. C 128), guanosine-5'-O-diphosphate (GDP-Gel, Cat. No. G 028) and uridine-5'-O-diphosphate (UDP-Gel, Cat. No. U 017).

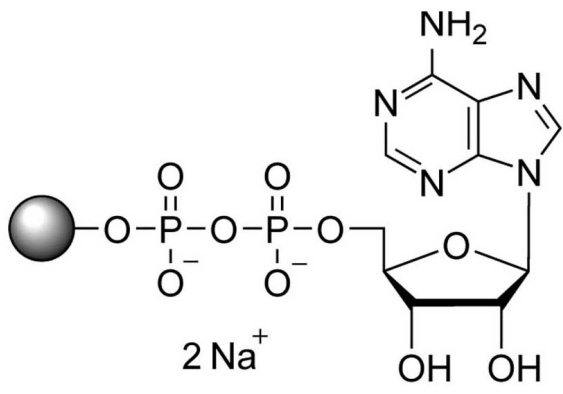
Ordering Information

[Order Online »](#)

BLG-A236-06	0.6ml
BLG-A236-25	2.5ml
BLG-A236-60	6ml

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO -Research Use Only

Product Details

Alternative Name Adenosine-5'-O-diphosphate, immobilized on a polymeric matrix

Couple Type Modified nucleotides

Formulation Suspension in 30 mM Na₂HPO₄ buffer (pH 7).

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com